

Product datasheet

Mouse Axl ELISA Kit ab207619

Recombinant SimpleStep ELISA

5 Images

Overview

**Product name** Mouse Axl ELISA Kit

**Detection method** Colorimetric

**Precision**

Intra-assay

Sample	n	Mean	SD	CV%
Serum	5			2.4%

Inter-assay

Sample	n	Mean	SD	CV%
Serum	3			3.4%

**Sample type** Cell culture supernatant, Urine, Serum, Hep Plasma, EDTA Plasma, Cit plasma

**Assay type** Sandwich (quantitative)

**Sensitivity** 10.8 pg/ml

**Range** 62.5 pg/ml - 4000 pg/ml

**Recovery**

Sample specific recovery

Sample type	Average %	Range
Cell culture supernatant	106	101% - 114%
Urine	109	104% - 114%
Serum	104	91% - 116%
Hep Plasma	83	74% - 91%
EDTA Plasma	100	86% - 111%
Cit plasma	102	90% - 111%

**Assay time** 1h 30m

**Assay duration** One step assay

**Species reactivity** **Reacts with:** Mouse  
**Does not react with:** Cow, Pig

**Product overview** Mouse Axl ELISA Kit (ab207619) is a single-wash 90 min sandwich ELISA designed for the quantitative measurement of Axl protein in cell culture supernatant, serum, urine, cit plasma, edta plasma, and hep plasma. It uses our proprietary SimpleStep ELISA® technology. Quantitate Mouse Axl with 10.8 pg/ml sensitivity.

SimpleStep ELISA® technology employs capture antibodies conjugated to an affinity tag that is recognized by the monoclonal antibody used to coat our SimpleStep ELISA® plates. This approach to sandwich ELISA allows the formation of the antibody-analyte sandwich complex in a single step, significantly reducing assay time. See the SimpleStep ELISA® protocol summary in the image section for further details. Our SimpleStep ELISA® technology provides several benefits:

- Single-wash protocol reduces assay time to 90 minutes or less
- High sensitivity, specificity and reproducibility from superior antibodies
- Fully validated in biological samples
- 96-wells plate breakable into 12 x 8 wells strips

A 384-well SimpleStep ELISA® microplate ([ab203359](#)) is available to use as an alternative to the 96-well microplate provided with SimpleStep ELISA® kits.

**Notes** Axl is a transmembrane receptor tyrosine kinase on the plasma membrane. The growth factor GAS6 is a ligand for Axl and regulates many physiological processes including cell survival, cell proliferation, migration and differentiation. The antibodies used in this kit were raised against a portion of the extracellular domain of Axl.

**Platform** Microplate (12 x 8 well strips)

**Properties**

**Storage instructions** Store at +4°C. Please refer to protocols.

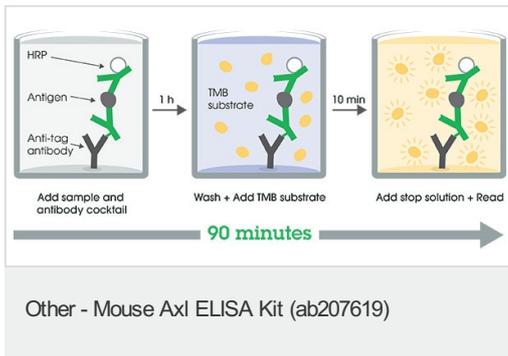
Components	1 x 96 tests
10X Mouse Axl Capture Antibody	1 x 600µl
10X Mouse Axl Detector Antibody	1 x 600µl
10X Wash Buffer PT (ab206977)	1 x 20ml
Antibody Diluent 5BR	1 x 6ml
Mouse Axl Lyophilized Recombinant Protein	2 vials
Plate Seals	1 unit

Components	1 x 96 tests
Sample Diluent NS (ab193972)	1 x 50ml
SimpleStep Pre-Coated 96-Well Microplate (ab206978)	1 unit
Stop Solution	1 x 12ml
TMB Development Solution	1 x 12ml

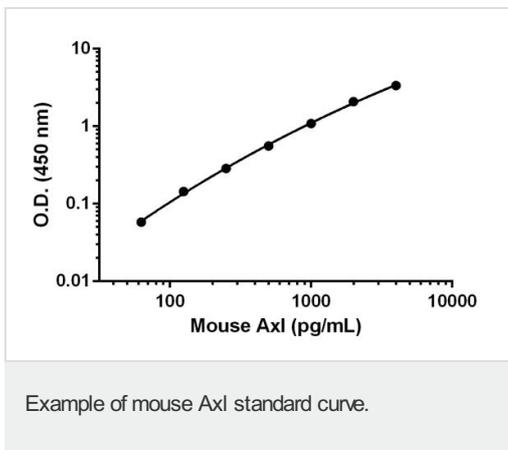
## Relevance

The protein encoded by this gene is a member of the receptor tyrosine kinase subfamily. Although it is similar to other receptor tyrosine kinases, the Axl protein represents a unique structure of the extracellular region that juxtaposes IgL and FNIII repeats. It transduces signals from the extracellular matrix into the cytoplasm by binding growth factors such as vitamin K dependent protein growth arrest specific gene 6. It is involved in the stimulation of cell proliferation. This receptor can also mediate cell aggregation by homophilic binding. Axl is a chronic myelogenous leukemia associated oncogene and also associated with colon cancer and melanoma. It is in close vicinity to the bcl3 oncogene which is at 19q13.1-q13.2. The Axl gene is evolutionarily conserved between vertebrate species. This gene has two different alternatively spliced transcript variants (AXL1 and AXL2).

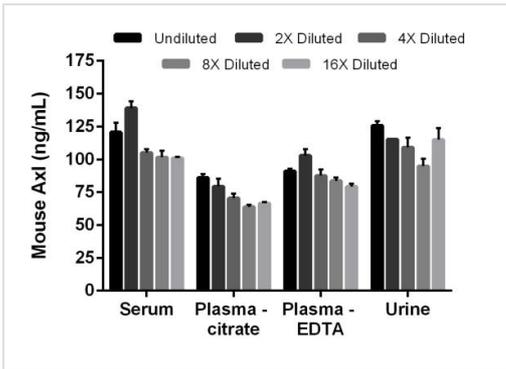
## Images



SimpleStep ELISA technology allows the formation of the antibody-antigen complex in one single step, reducing assay time to 90 minutes. Add samples or standards and antibody mix to wells all at once, incubate, wash, and add your final substrate. See protocol for a detailed step-by-step guide.

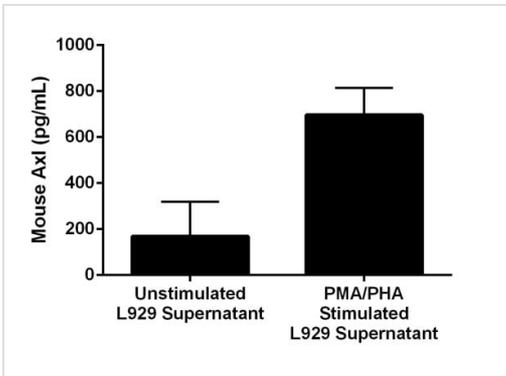


Background-subtracted data values (mean +/- SD) are graphed.



Interpolated concentrations of native Axl in mouse serum, plasma, and urine samples.

The concentrations of Axl were measured in duplicates, interpolated from the Axl standard curves and corrected for sample dilution. Undiluted samples are as follows: serum 4%, plasma (citrate) 4%, plasma (EDTA) 4%, and urine 1%. The interpolated dilution factor corrected values are plotted (mean +/- SD, n=2). The mean Axl concentration was determined to be 113 ng/mL in serum, 73.3 ng/mL in plasma (citrate), 88.9 ng/mL in plasma (heparin), and 112 ng/mL in urine.



Comparison of secreted Axl in unstimulated and PMA/PHA-stimulated L929 Cells.

L929 cells were grown in the absence (unstimulated) or presence of Phorbol Myristate Acetate (PMA) and phytohemagglutinin (PHA) (stimulated) for 3 days. Axl was measured in 2-fold diluted cell culture supernatants of unstimulated and PMA/PHA stimulated L929 and cell culture media. Measured values were interpolated from the Axl Standard Curve diluted in Sample Diluent NS and corrected for dilution factor. Mean of duplicate values +/-SD are graphed: 169 pg/mL unstimulated, 696 pg/mL stimulated, and undetectable in media.

Powered by recombinant antibodies

- Research with confidence: Consistent and reproducible results
- Long-term and scalable supply: Recombinant technology
- Success from the first experiment: Confirmed specificity
- Ethical standards compliant: Animal-free production

Sandwich ELISA - Mouse Axl ELISA Kit (ab207619)

To learn more about the advantages of recombinant antibodies see [here](#).

**Please note:** All products are "FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES"

## **Our Abpromise to you: Quality guaranteed and expert technical support**

---

- Replacement or refund for products not performing as stated on the datasheet
- Valid for 12 months from date of delivery
- Response to your inquiry within 24 hours
  
- We provide support in Chinese, English, French, German, Japanese and Spanish
- Extensive multi-media technical resources to help you
- We investigate all quality concerns to ensure our products perform to the highest standards

If the product does not perform as described on this datasheet, we will offer a refund or replacement. For full details of the Abpromise, please visit <https://www.abcam.com/abpromise> or contact our technical team.

## **Terms and conditions**

---

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors